

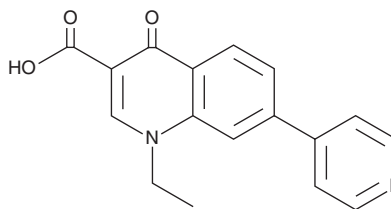
# PRODUCT INFORMATION



## Rosoxacin

Item No. 44519

**CAS Registry No.:** 40034-42-2  
**Formal Name:** 1-ethyl-1,4-dihydro-4-oxo-7-(4-pyridinyl)-3-quinolinecarboxylic acid  
**Synonym:** WIN 35,213  
**MF:** C<sub>17</sub>H<sub>14</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 294.3  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Rosoxacin is supplied as a solid. A stock solution may be made by dissolving the rososaxacin in the solvent of choice, which should be purged with an inert gas. Rosoxacin is slightly soluble (0.1-1 mg/ml) in acetonitrile.

### Description

Rosoxacin is a quinolone antibiotic.<sup>1</sup> It is active against *C. jejuni*, *S. boydii*, *S. dysenteriae*, *V. cholerae*, and *Y. enterocolitica* (MIC<sub>50</sub>s = ≤0.25 mg/L for all). Rosoxacin (10 μM) inhibits IL-1 receptor-associated kinase 1 (IRAK-1) *in vitro* and suppresses the development of murine growth hormone-induced experimental autoimmune hypophysitis (EAH) in mice when administered at a dose of 200 μg/animal per day.<sup>2</sup> Formulations containing rososaxacin have been used in the treatment of urinary tract infections and sexually transmitted diseases (STDs).

### References

1. Vanhoof, R., Hubrechts, J.M., Roebben, E., *et al.* The comparative activity of pefloxacin, enoxacin, ciprofloxacin and 13 other antimicrobial agents against enteropathogenic microorganisms. *Infection* **14**(6), 294-298 (1986).
2. Huang, H.C., Chen, Y.T., Lin, H.H., *et al.* Inhibition of IRAK1 is an effective therapy for autoimmune hypophysitis in mice. *Int. J. Mol. Sci.* **23**(23), 14958 (2022).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/17/2025

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
[WWW.CAYMANCHEM.COM](http://WWW.CAYMANCHEM.COM)